



EAGLE
BIOSCIENCES

Anti-Protein Kinase C Alpha α

Catalog Number: KCA00-A100

Aliquot: 100 μ L

Storage: -80°C

PRODUCT SPECIFICATION SHEET

- Type:** Polyclonal Rabbit IgG.
- Specificity:** Prepared against synthetic peptide corresponding to a distinct C terminal region, (SYVNPQFVHPILQSAV resides 657-672) from rat protein kinase C α , based on known homologies, is expected to react with most mammalian species.
- Applications:** Intended for use in Western blots, ELISA, Immunocytochemistry and Immunohistochemical techniques for fresh or frozen sections. Paraffin embedded and fixed tissues have not been tested.
- Recommended dilutions:** For Immunohistochemistry, Immunocytochemistry and Western blotting using alkaline phosphatase or horseradish peroxidase, a dilution of 1:250 is recommended. With a more sensitive detection system, such as enhanced chemiluminescence or ELISA, the antibody may be diluted further.
- Contents:** 100 μ L of rabbit antiserum containing 0.05% sodium azide.
- Storage:** Stable for several months at 4°C, up to a year at -20°C, and indefinitely at -80°C. Avoid repeated freeze thaws.
- References:**
1. Wetsel, W. C., et al., **1992** Tissue and Cellular Distribution of the Extended Family of Protein Kinase C Isoenzymes. *J. Cell Biol.* **117**: 21-133.
 2. *JBC* **268**:20110-20115.

Product Developed and Manufactured in the USA

EAGLE BIOSCIENCES, INC.
20A NW Blvd., Suite 112
Nashua, NH 03063

Phone: 617-419-2019
Fax: 617-419-1110
Email: info@EagleBio.com
www.EagleBio.com